



EXTENSION MEMO

Page 1 of 1

Product Description(s): S-PLEX® Human IL-10 Kit

Catalog Number(s) / Size(s): K151Y2S - Series

Lot Number: K00S0137

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K00S0137 has been extended to 31 DEC 2024.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
TURBO-BOOST™ Human IL-10 Antibody	D00T0033A	31 JUL 2024	31 DEC 2024
Biotin Human IL-10 Antibody	B00T0033A	31 DEC 2023	31 DEC 2024
Human IL-10 Calibrator	A00S0004	31 JUL 2024	31 DEC 2024
Diluent 2	M0020122	31 OCT 2024	31 DEC 2024
S-PLEX Coating Reagent C1	351987	30 APR 2024	31 DEC 2024
Blocker S1	R0010768	31 MAY 2024	31 DEC 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		20 OCT 2023

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: S-PLEX[®]Human IL-10 Kit

Catalog Number: K151Y2S - Series

Lot Number: K00S0137

Expiration Date: 31 Dec 2023

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
TURBO-BOOST [™] Human IL-10 Antibody	D00T0033A	2-8°C	31 Jul 2024
Biotin Human IL-10 Antibody	B00T0033A	2-8°C	31 Dec 2023
Human IL-10 Calibrator	A00S0004	2-8°C	31 Jul 2024
Diluent 2	M0020122	≤ -10°C	31 Oct 2024
Diluent 3	M0030117	≤ -10°C	31 Jan 2025
S-PLEX Coating Reagent C1	351987	≤ -70°C	30 Apr 2024
Blocker S1	R0010768	≤ -10°C	31 May 2024
S-PLEX Enhance E1	Y0550011	≤ -10°C	31 Jul 2025
S-PLEX Enhance E2	Y0560012	≤ -10°C	31 May 2025
S-PLEX Enhance E3	Y0570013	≤ -70°C	31 Jul 2025
S-PLEX Detect D1 with TURBO-TAG [™] reagent	Y0580016	≤ -70°C	31 May 2025
S-PLEX Detect D2	Y0590013	≤ -70°C	31 Jul 2025

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Name	Calibrator	
	Recovery	CV
IL-10	80% - 120%	< 20%

Name	Range Tested	LLOQ				LLOD	
		Observed	Units	Recovery	CV	Observed	Units
IL-10	2.45 - 39.2	9.80	fg/mL	80% - 120%	< 20%	1.23	fg/mL

Stock Concentration of Calibrator:

Name	Item	Concentration	Units
IL-10	20X Calibrator	380000	fg/mL

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		20 OCT 2023